

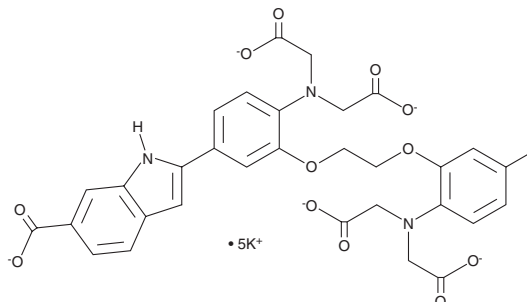
PRODUCT INFORMATION



Indo-1 (potassium salt)

Item No. 20418

CAS Registry No.: 132319-56-3
Formal Name: 2-[4-[bis(carboxymethyl)amino]-3-[2-[2-[bis(carboxymethyl)amino]-5-methylphenoxy]ethoxy]phenyl]-1H-indole-6-carboxylic acid, pentapotassium salt
MF: C₃₂H₂₆N₃O₁₂ • 5K
FW: 840.1
Purity: ≥90%
Ex./Em. Max: 349-364/475-485 (no Ca²⁺), 400-410 (Ca²⁺-bound) nm



Supplied as: A solid
Storage: -20°C
Stability: ≥2 years
Special Conditions: Protect from light

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Indo-1 (potassium salt) is supplied as a solid. It is soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Indo-1 is a ratiometric fluorescent calcium indicator. It is ideal for analyses using flow cytometry, as it uses a single excitation source, typically 349-364 nm light from an argon-ion laser. The emission maximum shifts from 475-485 nm without calcium to 400-410 nm when indo-1 binds calcium.^{1,2} Indo-1 is prone to photobleaching, which limits its usefulness in methods involving microscopy.¹

References

1. Paredes, R.M., Etzler, J.C., Watts, L.T., *et al.* Chemical calcium indicators. *Methods* **46(3)**, 143-151 (2008).
2. Grynkiewicz, G., Poenie, M., and Tsien, R.Y. A new generation of Ca²⁺ indicators with greatly improved fluorescence properties. *J. Biol. Chem.* **260(6)**, 3440-3450 (1985).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/11/2018

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM